

Product datasheet

Formaldehyde Assay Kit (Fluorometric) ab196997

[2 References](#) [2 Images](#)

Overview

Product name	Formaldehyde Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Saliva, Urine, Serum, Plasma, Other biological fluids, Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Sensitivity	> 2 μ M
Assay time	1h 00m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Formaldehyde Assay Kit (Fluorometric) (ab196997) provides a simple, sensitive and high-throughput adaptable assay that detects biologically relevant concentrations of formaldehyde (HCHO) in various fluids and tissues. The assay relies on the oxidation of formaldehyde to produce a stable fluorescent signal, which is directly proportional to the amount of formaldehyde found in samples.

The kit can detect less than 2 μ M of formaldehyde in a variety of samples.

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K805 PicoProbe™ Formaldehyde Fluorometric Assay Kit. K805-100 is the same size as the 100 test size of ab196997.</p> <p>Formaldehyde (HCHO) is the simplest aldehyde and one of the most often used organic compounds in industrial processes due to its high reactivity with a variety of chemicals. Endogenous formaldehyde can be produced in organisms via metabolic processes while exogenous formaldehyde can be absorbed after oral, dermal or inhalation exposure. formaldehyde is highly toxic due to its capacity to covalently bind to macromolecules, such as DNA, and is well known for its neurotoxicological effects. Exposure to relatively high levels of formaldehyde is believed to cause leukemia, nose and nasopharyngeal cancer, etc.; however the epidemiological evidence is unclear. It can also affect memory and learning capacity.</p>
--------------	---

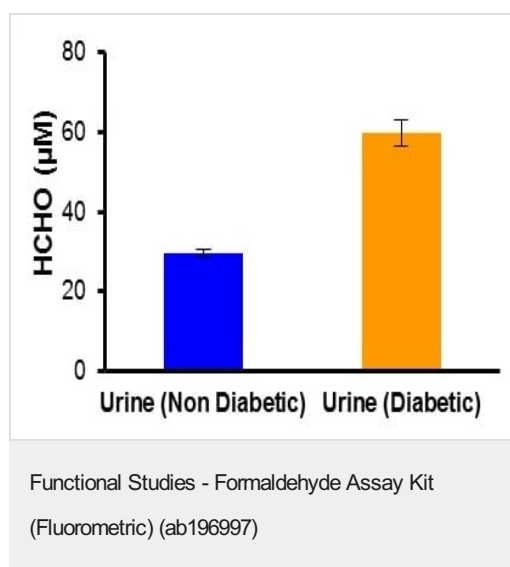
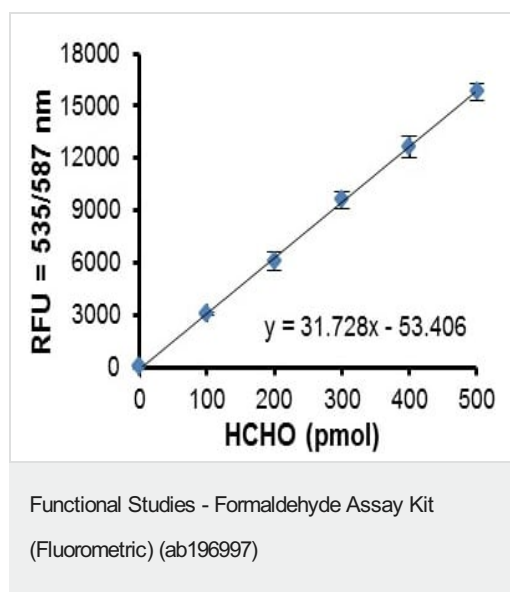
Platform	Microplate reader
-----------------	-------------------

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	100 tests
Developer Mix II	1 vial
Formaldehyde Enzyme Mix	1 vial
Formaldehyde Standard	1 x 100µl
HCHO Assay Buffer	1 x 25ml
PicoProbe I	1 x 0.4ml

Images



Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors